

Jumbo Star Sartopure® GF Plus

Adsorptive High Capacity
Filter Cartridge for Food &
Beverage Applications Only



Description

Jumbo Star Sartopure® GF Plus cartridges are optimized for the filtration of water and other beverages. Jumbo Star Sartoclean SC filter cartridges combine multiple layers of progressively finer glass fibre fleeces. Due to their high adsorptive power by glass fiber fleeces, they are an excellent solution for removal of particles and colloids from beverages.

Application

Typical applications for Jumbo Star Sartopure® GF Plus are:

- Prefiltration upstream of the final membrane filtration before filling
- Clarification step

Security

The Jumbo Star Sartopure® GF Plus filter elements ensure the selective and defined particle and colloid retention.

Performance

Jumbo Star Sartopure® GF Plus filter cartridges combine high dirt loading capacities with long service life and extremely high flow rates.

Flexibility

The modular concept allows the ideal adaptation to the customer requirements. Jumbo Star Sartopure® GF Plus filter cartridges are available in different heights.

Documentation

Jumbo Star Sartopure® GF filter cartridges are designed, developed and manufactured in accordance with a ISO 9001 Quality Management System. All materials used for the construction meet FDA requirements for use in food and beverage applications.

Specifications

Materials

Filter Material	Multiple glass fibre layers
Support Fleece	Polypropylene
Core	Polypropylene
End Caps	Polypropylene
O-Rings	Silicone

Retention Rates

0.65 µm
1.2 µm

Available Sizes | Filtration Area

Size 1	10"	5 m ²
Size 2	20"	10 m ²
Size 3	30"	15 m ²
Size 4	40"	20 m ²

Available Adapter

40

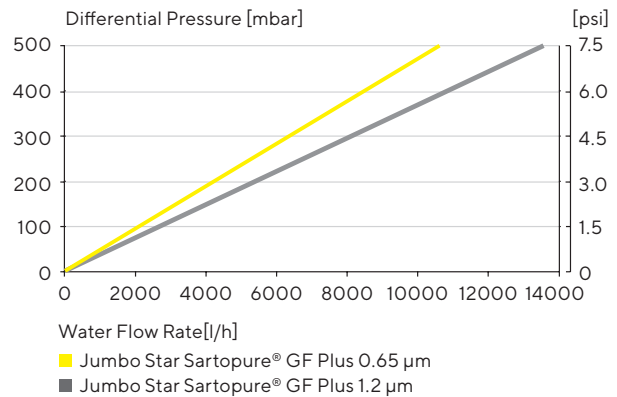
Operating Parameters

Max. Allowable Differential Pressure	4 bar 58 psi at 20 °C 1 bar 15 psi at 80 °C 0.5 bar 7.5 psi at 120 °C
Max. Allowable Back Pressure	1.5 bar 22 psi at 20 °C

Sterilisation

Hot Water	85-95 °C	30 min
In Line Steaming	110°C	30 min

Water Flow 10" Cartridge



Order Information

Order Code	Retention Rate [µm]
5554005J*	0.65 µm
5554003J*	1.20 µm

* = construction of height
1 = 250 mm (10")
2 = 500 mm (20")
3 = 750 mm (30")
4 = 1000 mm (40")

Germany

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Stedim North America Inc.
565 Johnson Avenue
Bohemia, NY 11716
Toll-Free +1 800 368 7178

For further contacts, visit
www.sartorius.com